

FMOC-Asp(EDANS)-OH *CAS 182253-73-2*

Catalog Number: 5000, 5001, 5002

Unit Size: 100 mg, 1 g, 5 g

Product Details

Storage Conditions Freeze (< -15 °C), Minimize light exposure

Expiration Date 12 months upon receiving

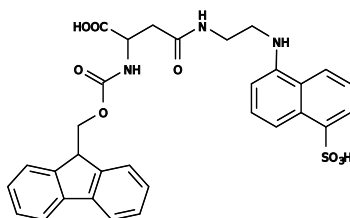
Chemical Properties

Appearance Solid grey

Molecular Weight 603.64

Soluble In DMF

Chemical Structure

**Spectral Properties**

Excitation Wavelength 336 nm

Emission Wavelength 455 nm

Applications

FMOC-Asp(EDANS)-OH is a building block for in-sequence peptide Asp labeling by EDANS. EDANS is one of the most common donor dyes for preparing FRET peptides (often paired with Dabcyl).